
argF Polyclonal Antibody

(Catalog #A54173)

Description

argF Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Escherichia coli

Isotype

IgG

Uniprot ID

P06960

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli Ornithine carbamoyltransferase subunit F protein (2-334AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

b0273, JW0266, argF, OTCas-2

Application

ELISA, Recommended dilution: ELISA 1:1000
